

PRODUCT INFORMATION



Pikromycin

Item No. 31331

CAS Registry No.: 19721-56-3
Formal Name: (3R,5R,6S,7S,9R,11E,13S,14R)-14-ethyl-13-hydroxy-3,5,7,9,13-pentamethyl-6-[[3,4,6-trideoxy-3-(dimethylamino)-β-D-xylohexopyranosyl]oxy]-oxacyclotetradec-11-ene-2,4,10-trione

Synonym: Picromycin

MF: C₂₈H₄₇NO₈

FW: 525.7

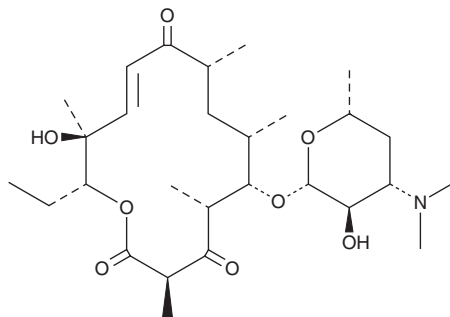
Purity: ≥95%

Supplied as: A crystalline solid

Storage: -20°C

Stability: ≥2 years

Item Origin: Bacterium/*Streptomyces* sp.



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Pikromycin is supplied as a crystalline solid. A stock solution may be made by dissolving the pikromycin in the solvent of choice, which should be purged with an inert gas. Pikromycin is soluble in DMSO.

Description

Pikromycin is a macrolide antibiotic that has been found in *S. venezuelae*.¹ It is active against *E. coli*, *S. aureus*, and *B. subtilis* (MIC_{99s} = 50-60, 90-100, and 25-30 μM, respectively).

Reference

1. Gupta, S., Lakshmanan, V., Kim, B.S., *et al.* Generation of novel pikromycin antibiotic products through mutasynthesis. *Chembiochem.* **9(10)**, 1609-1616 (2008).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 08/18/2020

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD

ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897

[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM

WWW.CAYMANCHEM.COM